



FGFR2

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : FGFR2

Substrate : TK

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[FGFR2] : 0.0123 ng/μL

Incubation time at RT : 50 min

ATP Km : 33.07 μM

Staurosporine IC₅₀ : 15.11 nM

